



PvuI (7)
SgfI (6) **MfeI (82)**
1 GGATCTGGATCGCTCCGGTGCCCGTCAGTGGGAGAGCGCACATCGCCACAGTCCCGGAGAAGTTGGGGGAGGGGTGGCAATTGAACGGGTGCCTA
101 GAGAAAGTGGCGCGGGTAAACTGGAAAGTGTGTCTGTACTGGCTCCGCCTTTTCCGAGGGTGGGGGAGAACCCTATATAAGTGCAGTAGTCGCC

HindIII (245)
Psp1406I (203) **PvuII (239)** **Bsu36I (291)**
201 GTGAACGTTCTTTTTCGCAACGGGTTTGCCGCCAGAACACAGCTGAAGCTTCGAGGGCTCGCATCTCTCTTACAGCGCCCGCCGCCCTACCTGAGGCC
301 GCCATCCACGCGGTTGAGTGCAGTCTGCCGCCTCCCGCCTGTGGTGCCTCTGAACTGCGTCCGCCGTCTAGGTAAGTTTAAAGCTCAGGTCGAGACC

NgoMIV (441)
401 GGGCCTTTGTCCGGCGCTCCCTTGAGCCTACCTAGACTCAGCGGCTCTCCACGCTTTGCTGACCCTGCTTGTCTAACTCTACGCTTTTGTTCGTTT

KasI (535) **AgeI (552)** **NcoI (560)**
BstEII (555)
501 TCTGTTCTGCGCCGTTACAGATCCAAGCTGTGACCGGGCGCTACCTGAGATCACCGGTCCACATGGACAAGTTTTGGTGGCACGCAGCCTGGGACTCTG
1▶ M D K F W W H A A W G L C

ClaI (625) **BbrPI (664)**
601 CCTCGTGGCGCTGAGCCTGGCGCAGATCGATTTGAATATAACCTGCCGCTTTCAGGTGTATTCCACGTGGAGAAAAATGGTGCCTACAGCATCTCTCGG
13▶ L V P L S L A Q I D L N I T C R F A G V F H V E K N G R Y S I S R

SdaI (788)
701 ACGGAGGCCGCTGACCTCTGCAAGCTTTCAATAGCACCTTGCCACAATGGCCAGATGGAGAAAGCTCTGAGCATCGGATTTGAGACCTGCAGGTATG
47▶ T E A A D L C K A F N S T L P T M A Q M E K A L S I G F E T C R Y

DraIII (814) **BamHI (830)**
BbrPI (814)
801 GGTTCATAGAAGGGCACGTGGTATCCCGGATCCACCCCACTCCATCTGTGCAGCAAACAACACAGGGGTGTACATCCTCACATCCAACACCTCCCA
80▶ G F I E G H V V I P R I H P N S I C A A N N T G V Y I L T S N T S Q
901 GTATGACACATATTGCTTCAATGCTTCAGCTCCACCTGAAGAAGATTGTACATCAGTCACAGACCTGCCAATGCCTTTGATGGACCAATTACCATACT
113▶ Y D T Y C F N A S A P P E E D C T S V T D L P N A F D G P I T I T

HpaI (1003)
1001 ATGTTAACCGTGATGGCACCCGCTATGTCCAGAAAGGAGAATACAGAACGAATCCTGAAGACATCTACCCAGCAACCCTACTGATGATGACGTGAGCA
147▶ I V N R D G T R Y V Q K G E Y R T N P E D I Y P S N P T D D D V S
1101 GCGGCTCTCCAGTGAAGGAGCAGCACTTCAGGAGTTACATCTTTTACACCTTTTCTACTGTACACCCATCCAGACGAAGACAGTCCCTGGATCAC
180▶ S G S S S E R S S T S G G Y I F Y T F S T V H P I P D E D S P W I T

XcmI (1263)
SmaI (1258)
BstXI (1245) **EcoO109I (1258)** **BsaBI (1277)**
1201 CGACAGCAGACAGAATCCCTGTACCA GAGACCAAGACACATTCCACCCAGTGGGGGTCCCATACCACTCATGGATCTGAATCAGATGGACTCA
213▶ D S T D R I P A T R D Q D T F H P S G G S H T T H G S E S D G H S
1301 CATGGGAGTCAAGAAGGTGGAGCAAACAACCTCTGGTCTATAAGGACACCCAAATCCAGAATGGCTGATCATCTTGGCATCCCTTTGGCCTTGG
247▶ H G S Q E G G A N T T S G P I R T P Q I P E W L I I L A S L L A L
1401 CTTTGATTCTGCAGTTTGCATTGCAGTCAACAGTCAAGAAGGTGTGGGAGAAAGAAAGCTAGTGATCAACAGTGGCAATGGAGCTGTGGAGGACAG
280▶ A L I L A V C I A V N S R R R C G Q K K K L V I N S G N G A V E D R

PvuII (1595)
1501 AAAGCCAAGTGGACTCAACGGAGAGGCCAGCAAGTCTCAGGAAATGGTGCATTTGGTGAACAAGGAGTCGTCAGAAACTCCAGACCAGTTTATGACAGCT
313▶ K P S G L N G E A S K S Q E M V H L V N K E S S E T P D Q F M T A

MscI (1656)
NheI (1650)
1601 GATGAGACAAGAACCTGCAGAATGTGGACATGAAGATTGGGGTGAACAGCTAGCTGGCCAGACATGATAAGATACATTGATGAGTTTGGACAAACCAC
347▶ D E T R N L Q N V D M K I G V •

HpaI (1788) **MfeI (1799)**
1701 AACTAGAATGCAGTGAATAAATGCTTTATTTGTGAAATTTGTGATGCTATTGCTTTATTTGTAACCATTATAAGCTGCAATAAACAAGTTAAACAACAAC

EcoRI (1884)
1801 AATTGCATTCATTTTATGTTTCAGGTTCCAGGGGAGGTGTGGGAGTTTTTAAAGCAAGTAAACCTCTACAAATGTGGTATGGAATTCTAAAATACAG
1901 CATAGCAAACTTTAACCTCAAATCAAGCCTCTACTTGAATCCTTTTCTGAGGGATGAATAAGGCATAGGCATCAGGGGCTGTTGCCAATGTGCATTAG

SapI (2066)
2001 CTGTTTGCAGCCTCACCTCTTTTCATGGAGTTTAAAGATATAGTGTATTTTCCCAAGTTTGAAGTCTCTTCTTTCTTTATGTTTTAAATGCACTGAC

SspI (2123) **SwaI (2137)** **EcoO109I**
2101 CTCCACATTCCTTTTTAGTAAATATTGAGAATAATTTAAATACATCATTGCAATGAAAATAAATGTTTTTATTAGGCAGAATCCAGATGCTCAAG
2201 GCCCTTCATAATATCCCCAGTTTAGTAGTTGGACTTAGGGAACAAGGAACCTTAAATAGAATTGGACAGCAAGAAAGCGAGCTTCTAGCTTTAGTTC
141▶ • N

SacI (2398)
2301 CTGGTGTACTTGAGGGGATGAGTTCCTCAATGGTGGTTTTGACCAGCTTGCCATTCTCTCAATGAGCACAAGCAGTCAGGAGCATAGTCAGAGATGA
138▶ R T Y K L P I L E E I T T K V L K G N M E I L V F C D P A Y D S I L

BstXI (2427)
2401 GCTCTCTGCACATGCCACAGGGGCTGACCACCTGATGGATCTGTCCACCTCATCAGAGTAGGGGTGCCTGACAGCCACAATGGTGTCAAAGTCTTCTG
105▶ E R C M G C P S V V R I S R D V E D S Y P H R V A V I T D F D K Q

